

## Rabbit Anti-Alexa Fluor 750 Antibody, Biotinylated

Catalog Number: 46610

Unit Size: 100 ug

### Product Details

---

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	1 mg/mL
Formulation	PBS, 0.2% (w/v) BSA

### Unit Details

---

Reconstitution	0.1 mL ddH <sub>2</sub> O
Volume	

### Antibody Properties

---

Clonality	Polyclonal
Host	Rabbit
Immunogen	Alexa Fluor 750 coupled to carrier protein

### Biological Properties

---

Stabilizer	BSA
Appearance	Solid
Preparation	It is produced with Alexa Fluor 750 coupled carrier protein in rabbit and affinity purified with Protein G beads. The antibody is conjugated with Biotin under optimal conditions.
Soluble In	Water
Application	IF, ELISA, Flow Cytometry (FACS)

### Spectral Properties

---

Conjugate	Biotin
-----------	--------

### Applications

---

Rabbit Anti-Alexa Fluor 750 Antibody, Biotinylated is a polyclonal antibody that specifically binds to Alexa Fluor® 750-conjugated molecules and is directly labeled with biotin. This format enables flexible detection using streptavidin- or avidin-based systems across a wide range of assay types.

The antibody may cross-react with related near-infrared fluorophores such as iFluor® 750, ATTO 740, and DyLight® 755. Its broad compatibility makes it ideal for signal amplification, conjugation verification, and multiplexed fluorescence detection making it a versatile tool for imaging and analytical workflows.

*Alexa Fluor® and DyLight® are registered trademarks of Thermo Fisher Scientific.*